

CNGB1 Antibody [#C15265]

Catalog Number:	C15265
Concentration:	1mg/ml
Swiss-Prot No.:	Q14028
Other Names:	CNCG4; CNG channel 4; CNG-4; CNG4; Cyclic nucleotide-gated cation channel modulatory subunit
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human CNGB1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CNGB1 antibody detects endogenous levels of total CNGB1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%) Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Chen T.Y., Nature 362:764-767(1993). Bradley J., Nat. Neurosci. 7:705-710(2004).